

Batch information

Diluent 5 HTRF Diluent 5 (5X) - 2 mL
Product reference : **62DL5DDA**

Product characteristics :

- Lot Number 16G
- Conditioning buffer -
- Appearance Frozen form
- Storage conditions <-16 °C
- Expiration date 2025.06.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	2.00 mL	N/A

PLEASE NOTE

The expiration date fo batch 16G has been extended to 2025-06-19 after validation by QC.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

